

PRODUCT INFORMATION



(5Z,2E)-CU-3

Item No. 37770

CAS Registry No.: 1815598-71-0
Formal Name: N-[(5Z)-5-[(2E)-3-(2-furanyl)-2-propen-1-ylidene]-4-oxo-2-thioxo-3-thiazolidinyl]-benzenesulfonamide

MF: C₁₆H₁₂N₂O₄S₃

FW: 392.5

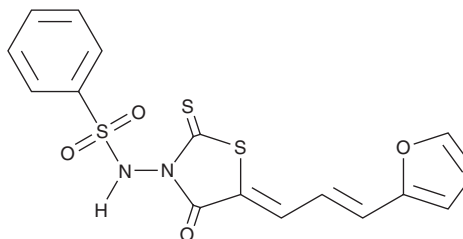
Purity: ≥98%

UV/Vis.: λ_{max}: 419 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

(5Z,2E)-CU-3 is supplied as a solid. A stock solution may be made by dissolving the (5Z,2E)-CU-3 in the solvent of choice, which should be purged with an inert gas. (5Z,2E)-CU-3 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of (5Z,2E)-CU-3 in these solvents is approximately 5 and 2 mg/ml, respectively.

Description

(5Z,2E)-CU-3 is an inhibitor of diacylglycerol kinase α (DGK- α ; IC₅₀ = 0.6 μ M for the porcine enzyme).¹ It is selective for DGK- α over other type I DGKs, as well as type II, type III, type IV, and type V DGKs (IC₅₀s = 7-36 μ M). (5Z,2E)-CU-3 (5 μ M) induces apoptosis in HeLa cervical and HepG2 liver cancer cells and enhances increases in *IL2* mRNA expression in Jurkat T cells induced with concanavalin A (Item No. 14951).

Reference

1. Liu, K., Kunii, N., Sakuma, M., *et al.* A novel diacylglycerol kinase α -selective inhibitor, CU-3, induces cancer cell apoptosis and enhances immune response. *J. Lipid Res.* **57(3)**, 368-379 (2016).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/25/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM